

Nitric oxide synthase (endothelial) polyclonal antibody

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-505-R100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name eNOS, NOS III

Application ICC, WB

Application Notes Immunohistochemical staining of eNOS (NOS III) in mouse cardiac tissue with Prod. No. ALX-210-505 results in the staining of the endothelium.
Detects a band of ~140kDa by Western blot.

Crossreactivity Does not cross-react with inducible NOS (iNOS; NOS II) or neuronal NOS (nNOS; NOS I).

Formulation Liquid. Neat serum diluted in PBS containing 0.05% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 598-612 (P598YNSSPRPEQHKSYK612C) of bovine endothelial nitric oxide synthase (eNOS; NOS III).

Recommendation Immunocytochemistry (1:250)Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Dog, Human, Mouse, Rat

UniProt ID P29473

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)